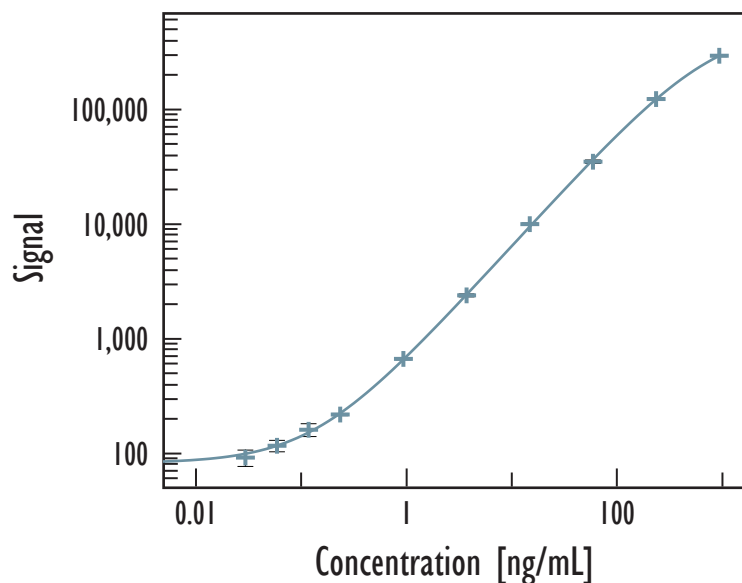


# MULTI-ARRAY<sup>®</sup> Thrombomodulin Assay

## Detection of Thrombomodulin in Serum and Plasma Samples



Standard curve data is from a representative experiment

Detection Limits (ng/mL)	Thrombomodulin
Average LLOD	0.08
Average LLOQ	0.8

LLOD is defined as 2.5x stdev above the background

LLOQ is designated by the lowest concentration on the standard curve where % CV is less than 20% and recovery of predicted concentration is within 20% of 100%

Kit Size	Catalog Number
1 plate	K151EKC-1
5 plates	K151EKC-2
20 plates	K151EKC-3
20 plates (Base)	K151EKA-3

### Endogenous Levels in Human Samples

- 16 normal human donors; matched sera and plasmas
- Detected level was above LLOQ in all samples
- Average CVs for measured samples was less than 10% for all assays

		ng/mL
Serum	Mean	2.2
	Median	2.2
	Range	1.3 - 3.0
EDTA Plasma	Mean	1.4
	Median	1.3
	Range	0.9 - 2.1
Heparin Plasma	Mean	1.9
	Median	1.9
	Range	1.2 - 3.6

# MULTI-ARRAY<sup>®</sup> Thrombomodulin Assay

## Detection of Thrombomodulin in Serum and Plasma Samples

### Dilutional Linearity

- Measured endogenous analyte levels in samples diluted into assay diluent

$$\% \text{ recovery} = \frac{(\text{measured value} * \text{dilution factor} * 100)}{\text{predicted value}}$$

	Dilution Factor	% Recovery
Serum	1/2	107
	1/4	99
	1/8	110
EDTA Plasma	1/2	120
	1/4	111
	1/8	114
Heparin Plasma	1/2	114
	1/4	108
	1/8	105

Values presented are averages across two pooled samples

### Spike Recovery

- Measured analyte spiked into undiluted human samples

$$\% \text{ recovery} = \frac{(\text{measured value} * 100)}{\text{expected value}}$$

	% Recovery
Serum	87
EDTA Plasma	69
Heparin Plasma	83

Values presented are averages across two pooled samples and including spike levels of 12.5, 25, and 50 ng/mL